

RET (M918T) Kinase Assay Service

Activity-based RET (M918T) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

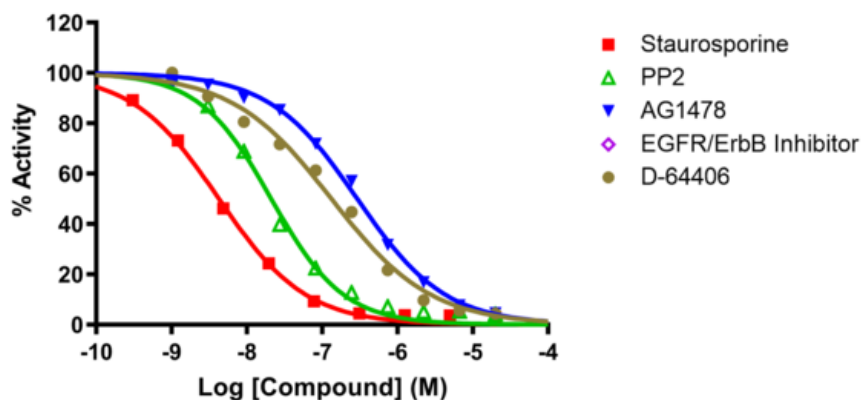
Target Overview

| | |
|-------------------|--|
| CAT# | RET (M918T) |
| Enzyme | Human RET (M918T) |
| HGNC | RET (M918T) |
| Synonym | CDHF12, CDHR16, HSCR1, MEN2A, MEN2B, MTC1, PTC, RET51 |
| Substrates | Peptide substrate, [RRRRRRRRRRRVYSTDYRLFNPS], 20 μM ATP 10 μM |

Assay Properties

| | | |
|---------------------------------|--|------------------|
| Reaction | Substrate + [³² P]-ATP = ³² P-Substrate + ADP | |
| Readout | Enzymatic activity | |
| Reference Compound IC50s | Compounds | IC50 (nM) |
| | Staurosporine | 4.3 |
| | PP2 | 21 |
| | AG1478 | 290 |
| | EGFR/ErbB Inhibitor | >100,000 |
| | D-64406 | 130 |
| Scientific Information | P07949 (http://www.uniprot.org/uniprot/P07949) | |
| Screening Location | Malvern, USA | |

Reference compound IC50 for RET (M918T)



Let's Discover Together.

© Aug 2022 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/ret_m918t_kin_malvern

